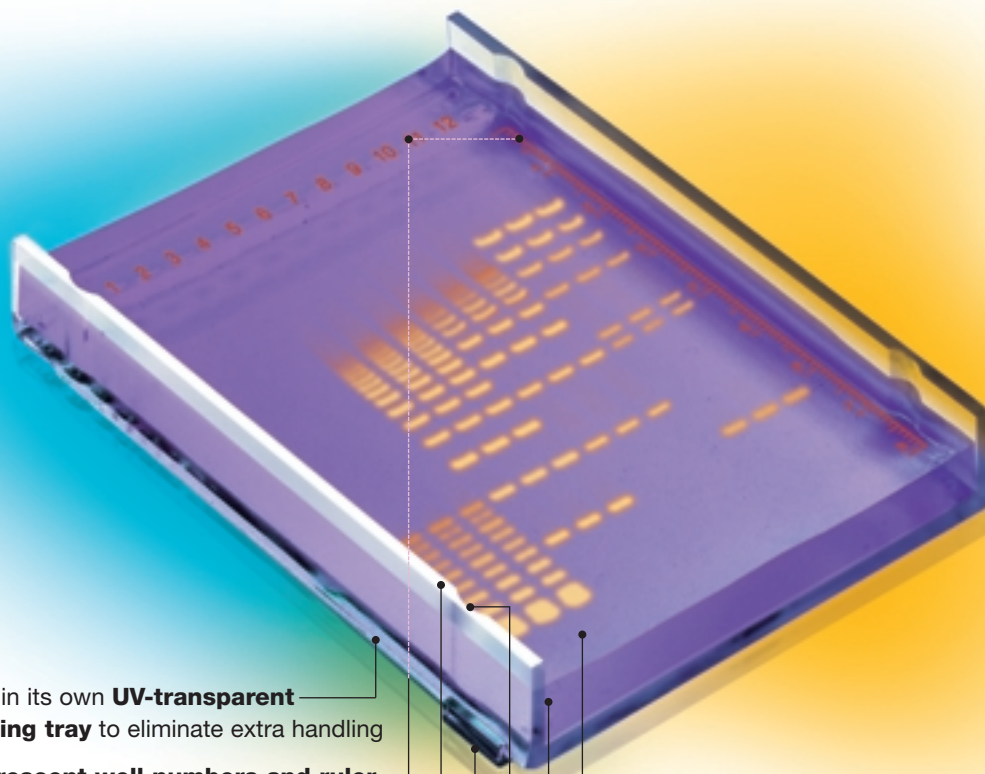


Get Unrivaled Performance

with **ReadyAgarose™** Gels



Cast in its own **UV-transparent running tray** to eliminate extra handling

Fluorescent well numbers and ruler printed on tray for easy reference

Two-sided tray for consistently straight lanes and familiar run conditions

Lock-tabs on patent-pending tray prevent tray movement during loading and running

Supplied in **TBE or TAE buffer**, with or without ethidium bromide — gels that fit your needs

Offered in **mini and wide mini gel formats** and 8, 12, 20, 32, or 64 wells for all loading volume needs

Cast with **high-strength, high-resolution agaroses** that are qualified for molecular biology applications

Available in **0.8%, 1.0%, and 3.0% agarose** to separate DNA from 20 to 20,000 bp

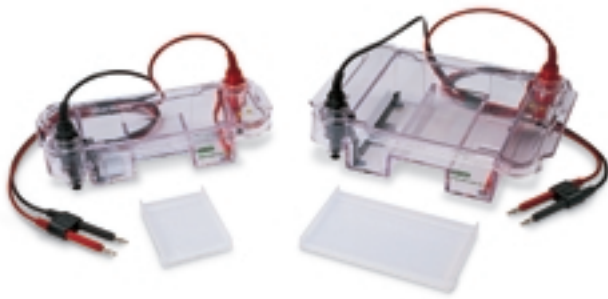
Notched tray holds gel in place

Use the gel width ruler below to measure the inside dimension of your subcell. Mini gels fit every subcell at least 7.1 cm wide; wide mini gels fit every subcell at least 15.6 cm wide

BIO-RAD

mini gel tray width

wide mini gel tray width



ReadyAgarose Gels

You want unequaled performance —

The new ReadyAgarose precast gels are the answer. The UV-transparent gel tray allows you to see and record results without handling the gel. The 2-sided tray design ensures straight and reproducible lanes. Also, ReadyAgarose gels are made with Bio-Rad's high-strength, high-resolution Certified™ agaroses to meet the needs of your molecular biology applications.

You want easy handling and tracking —

Bio-Rad's unique lock-tab design secures the tray in Sub-Cell® GT cells during loading and running. Fluorescent well numbers and ruler are printed on the tray for easy loading and tracking of results.

You want gels to accommodate all your sample and subcell requirements —

Our 57 gel types give you a wide choice of well numbers and sample capacity. ReadyAgarose gels are designed to fit almost every subcell available. Use the gel width ruler on the front page to see whether ReadyAgarose gels fit your system.

Bio-Rad's ReadyAgarose precast gels provide unrivaled performance — just what you need for superior results. To request a free sample, contact your local Bio-Rad representative or visit us on the Web at www.bio-rad.com/readyagarose/

Ordering Information

	ReadyAgarose Mini Gels			ReadyAgarose Wide Mini Gels		
	8-well (20 µl)	12-well (10 µl)	2 x 8-well (20 µl)	20-well (20 µl)	32-well* (15 µl)	2 x 32-well* (15 µl)
TBE Gels						
0.8%	161-3001	161-3007	—	161-3025	161-3031	—
0.8% plus ethidium bromide	161-3002	161-3008	—	161-3026	161-3032	—
1%	161-3003	161-3009	—	161-3027	161-3033	161-3037
1% plus ethidium bromide	161-3004	161-3010	—	161-3028	161-3034	161-3038
3%	161-3005	161-3011	—	161-3029	161-3035	161-3039
3% plus ethidium bromide	161-3006	161-3012	—	161-3030	161-3036	161-3040
TAE Gels						
0.8%	161-3013	161-3019	—	161-3041	161-3047	—
0.8% plus ethidium bromide	161-3014	161-3020	—	161-3042	161-3048	—
1%	161-3015	161-3021	161-3057	161-3043	161-3049	161-3053
1% plus ethidium bromide	161-3016	161-3022	—	161-3044	161-3050	161-3054
3%	161-3017	161-3023	—	161-3045	161-3051	161-3055
3% plus ethidium bromide	161-3018	161-3024	—	161-3046	161-3052	161-3056

Gel shelf life at room temperature: 12 months without ethidium bromide; 6 months with ethidium bromide.
*Accommodates multichannel pipets.

Catalog # Description

Molecular Rulers

170-8351	20 bp EZ Load™ Molecular Ruler, 20–1,000 bp range, 100 applications
170-8352	100 bp EZ Load Molecular Ruler, 100–1,000 bp range, 100 applications
170-8200	AmpliSize® Molecular Ruler, 50–2,000 bp range, 50 applications
170-8353	100 bp PCR EZ Load Molecular Ruler, 100–3,000 bp range, 100 applications
170-8354	500 bp EZ Load Molecular Ruler, 500–8,000 bp range, 100 applications
170-8355	1 kb EZ Load Molecular Ruler, 1–15 kb range, 100 applications
170-8205	2.5 kb Molecular Ruler, 2.5–35 kb range, 100 applications
170-8356	EZ Load Precision Molecular Mass Ruler, 100–1,000 bp, 10–100 ng, 100 applications

Premixed Buffers

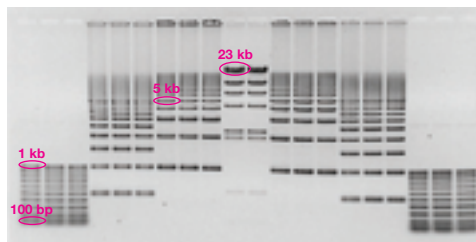
161-0767	5x Nucleic Acid Sample Buffer, 10 ml
161-0733	10x TBE, 1 L
161-0770	10x TBE, 5 L cube
161-0743	50x TAE, 1 L
161-0773	50x TAE, 5 L cube

Ethidium Bromide Solutions and Tablets

161-0433	Ethidium Bromide Solution, 10 ml, 10 mg/ml
161-0430	Ethidium Bromide Tablets, 10 x 11 mg

ReadySub-Cell™ Electrophoresis Cells

170-4487	Mini ReadySub-Cell GT Cell, for ReadyAgarose gels, includes subcell unit only
170-4489	Wide Mini ReadySub-Cell GT Cell, for ReadyAgarose gels, includes subcell unit only



ReadyAgarose 1% 20-well TAE gel.

Repeated loading of 100 bp, 500 bp, and 1 kb molecular rulers, and lambda HindIII digest. Run at 75 V for 80 min with 1x TAE running buffer.

	Mini	Wide Mini
Tray dimensions	7.2 x 10 cm	15.6 x 10 cm
Gel dimensions	6.6 x 10 cm	15.1 x 10 cm

Gel Selection Guide

Gel	Resolution Range
0.8%	300 bp–20 kb
1.0%	200 bp–10 kb
3.0%**	20 bp–1,000 bp

**Our 3% gel is made with high-resolution agarose, which makes it easier to handle and provides the same resolution as a standard 4% agarose gel.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191
Canada (905) 712-2771 China (86-21) 63052255 Denmark 45 44 52-1000 Finland 358 (0)9 804 2200 France 01 47 95 69 65 Germany 089 318 84-177
Hong Kong 852-2789-3300 India (91-124) 6398112/113/114, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270
Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64-9-4152280
Norway 47-23-38-41-30 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-2729877 South Africa 00 27 11 4428508 Spain 590 5200
Sweden 46 (0)8-55 51 27 00 Switzerland 061 717-9555 United Kingdom 0800-181134